

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3419P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the carboxy terminus of Anthrax edema factor
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
<b>Storage &amp; Stability</b>	Store Anthrax Edema Factor peptide at -20°C, stable for one year.
<b>Application</b>	Anthrax peptide is used for blocking the activity of Anthrax edema factor antibody.
<b>References</b>	<p>Robertson DL, Tippetts MT, and Leppla SH. Nucleotide sequence of the Bacillus anthracis edema factor gene (cya): a calmodulin-dependent adenylate cyclase. <i>Gene</i> 1988; 73:363-71.</p> <p>Lanahan A, Williams JB, Sanders LK, et al. Growth factor-induced delayed early response genes. <i>Mol Cell Biol.</i> 1992; 12:3919-29.</p> <p>Sweet MJ, Leung BP, Kang D, et al. A novel pathway regulating lipopolysaccharide-induced shock by ST2/T1 via inhibition of Toll-like receptor 4 expression. <i>J. Immunol.</i> 2001; 166:6633-9.</p> <p>Wald D, Qin J, Zhao Z, et al. SIGIRR, a negative regulator of Toll-like receptor-interleukin 1 receptor signaling. <i>Nat. Immunol.</i> 2003; 4:920-7.</p>

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.